

PRODUCT INFORMATION



17(R)-Protectin D1

Item No. 34198

CAS Registry No.: 1365694-03-6
Formal Name: 10R,17R-dihydroxy-4Z,7Z,11E,13E,15Z,19Z-docosahexaenoic acid

Synonyms: Aspirin-triggered Protectin D1, 17-*epi* Neuroprotectin D1, 17(R)-Neuroprotectin D1, 17(R)-NPD1, 17(R)-PD1, 17-*epi* Protectin D1

MF: C₂₂H₃₂O₄

FW: 360.5

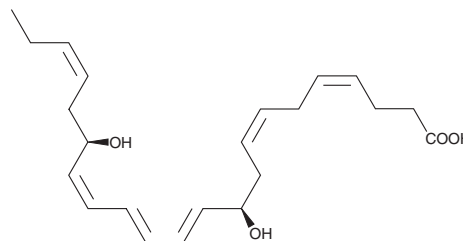
Purity: ≥98%

UV/Vis.: λ_{max}: 272, 283 nm

Supplied as: A solution in ethanol

Storage: -80°C

Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

17(R)-Protectin D1 is an aspirin-triggered epimer of the specialized pro-resolving mediator protectin D1 (Item No. 10010390).¹ It decreases leukotriene B₄-induced transendothelial migration of human polymorphonuclear (PMN) neutrophils when used at concentrations ranging from 0.1 to 10 nM. 17(R)-Protectin D1 (0.01-10 ng) reduces the recruitment of neutrophils in a mouse model of TNF-α-induced peritonitis.

Reference

1. Serhan, C.N., Fredman, G., Yang, R., *et al.* Novel proresolving aspirin-triggered DHA pathway. *Chem. Biol.* **18**(18), 976-987 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/21/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM